

SFX succinimidyl ester, 5 isomer

5-SFX has a 6-atom aminohexanoyl spacer group that separates the fluorescein dye and the amine-reactive group.



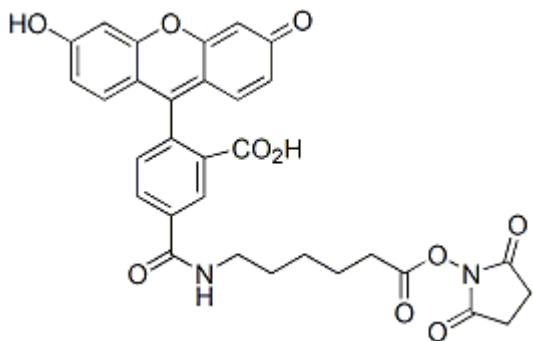
Product attributes

Excitation/Emission	494/521 nm (pH 9)
---------------------	-------------------

Product Description

5-SFX (full name: 6-(Fluorescein-5-carboxamido)hexanoic acid, succinimidyl ester, single isomer) has a 6-atom aminohexanoyl spacer group that separates the fluorescein dye and the amine-reactive group.

- $\lambda_{Ex}/\lambda_{Em}$ (pH 9) = 494/521 nm
- Yellow solid soluble in DMF, DMSO or pH>6.5 buffers
- Store at -20 °C and protect from light
- $C_{31}H_{26}N_2O_{10}$
- MW:586.55



This datasheet was generated on December 16, 2025 at 12:06:24 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5-sfx-6-fluorescein-5-carboxamidohexanoic-acid-succinimidyl-estersingle-isomer/>